

PowerPac Adapter Catalog number 165-5061



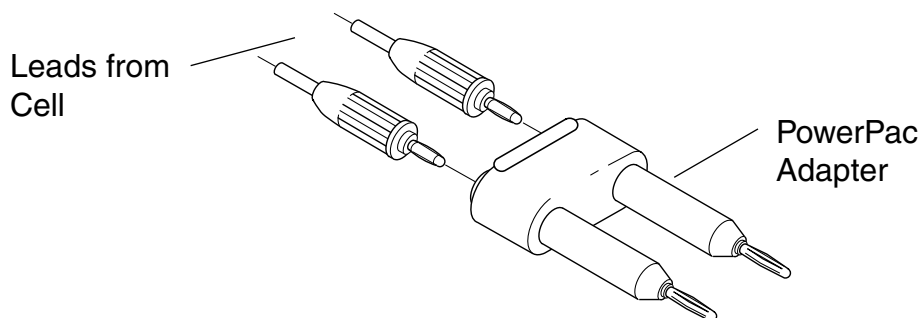
Bio-Rad PowerPac power supplies are designed and certified to conform to IEC1010-1, an internationally accepted standard for electrical safety of laboratory instrumentation. Certified products are safe to use when operated in accordance with the instruction manual provided with the product. The safety certification does not extend to electrophoresis cells or accessories, which are not IEC1010-1 certified, even when connected to the power supply.

This PowerPac Adapter is **NOT** certified to conform with IEC1010-1. Some non-certified electrophoresis cells, with leads less than 17 mm in length, may require the use of this adapter. Use of this adapter with any Bio-Rad PowerPac power supply will negate the IEC1010-1 safety certification, and may present a safety hazard to the user.

Instructions:

The following guidelines should be observed and followed when using the Bio-Rad PowerPac Adapter:

- Keep the PowerPac Adapter connected to the electrophoresis cell leads rather than the PowerPac power supply.
- Follow all safety precautions as listed in the safety section of the PowerPac instruction manual.



1. Insert the leads for the non-certified electrophoresis cells into the PowerPac Adapter. Ensure that the polarity of the cables is maintained (*i.e.*, red to red and black to black).
2. Insert the PowerPac Adapter into the correct output terminals of the PowerPac power supply maintaining the polarity of the cables (*i.e.*, red to red and black to black).